
BHMT2 Polyclonal Antibody

(Catalog #A64140)

Background

Involved in the regulation of homocysteine metabolism. Converts homocysteine to methionine using S-methylmethionine (SMM) as a methyl donor.

Description

BHMT2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9H2M3

Purification

>95%, Protein G purified

Immunogen

Recombinant Human S-methylmethionine--homocysteine S-methyltransferase BHMT2 protein (1-200AA)

Storage

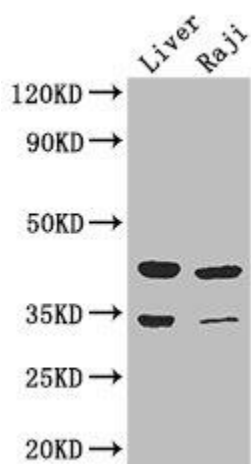
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

S-methylmethionine--homocysteine S-methyltransferase BHMT2, SMM-hcy methyltransferase, Betaine--homocysteine S-methyltransferase 2, BHMT2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

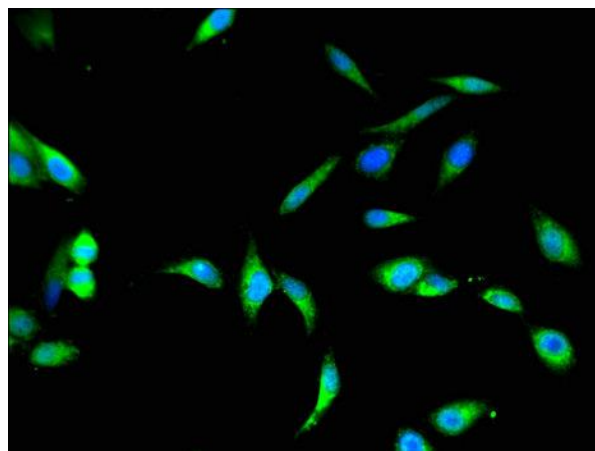
Positive WB detected in: Mouse liver tissue, Raji whole cell lysate

All lanes: BHMT2 Polyclonal Antibody at 3ug/ml
Secondary

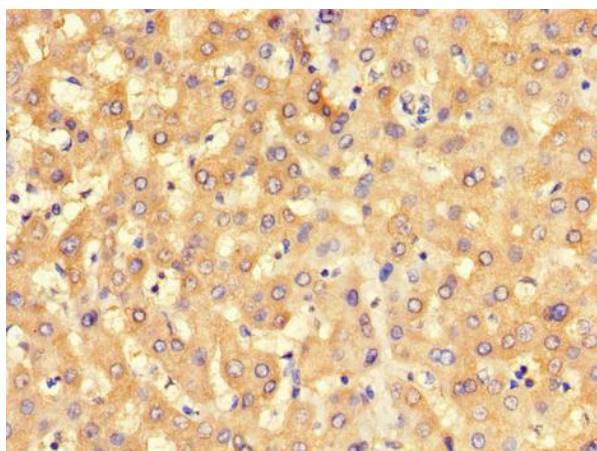
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 41, 34 kDa

Observed band size: 41, 34 kDa



Immunofluorescent analysis of HeLa cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human liver tissue at dilution of 1:100